

Biotin-4-fluorescein

Biotin-4-fluorescein (*N*-(5-aminopentyl)biotinamide, trifluoroacetic acid salt) has been reported to be a superior fluorescent biotin derivative for accurate measurement of avidin and streptavidin in crude biological samples. Binding of biotin to avidin and streptavidin is high affinity, fast and noncooperative.



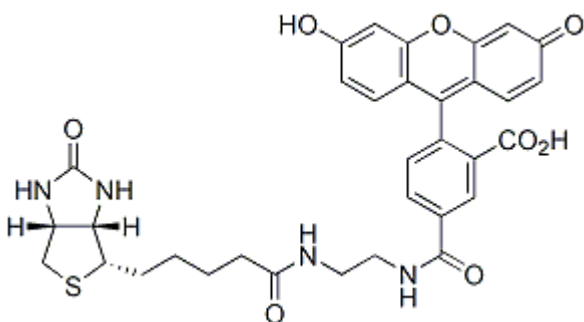
Product attributes

Excitation/Emission	494/523 nm
CAS number	1032732-74-3

Product Description

Biotin-4-fluorescein (*N*-(5-aminopentyl)biotinamide, trifluoroacetic acid salt) has been reported to be a superior fluorescent biotin derivative for accurate measurement of avidin and streptavidin in crude biological samples. Binding of biotin to avidin and streptavidin is high affinity, fast and noncooperative.

- $\lambda_{Ex}/\lambda_{Em} = 494/523$ nm
- Yellow solid soluble in DMSO, or pH>7 buffers
- Store at -10 °C to -35 °C and protect from light
- $C_{33}H_{32}N_4O_8S$
- MW: 644.7



References

1. Biochim Biophys Acta 1427, 33 (1999).
2. Biochim Biophys Acta 1427, 44 (1999).

This datasheet was generated on June 5, 2026 at 07:19:35 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/biotin-4-fluorescein/>